

HTRF® Research Reagents

HTRF Human VHL Bind. Kit - 10K pts

Product Number : 64BDVHLPEH Lot Number : 02C

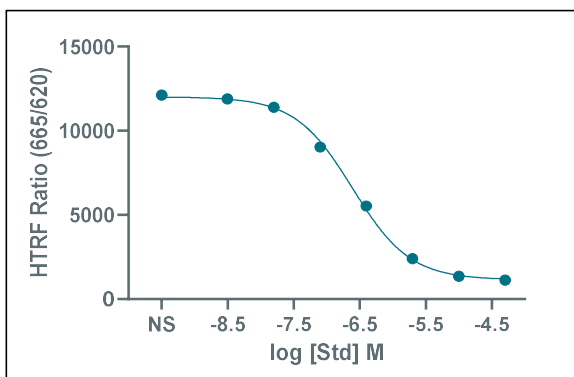
Expiration date : 2026-01-04 Document version : 1

Reagents :	Lot n°
HTRF Diluent 9 (5X) - 100 mL	04RA
HTRF PROTAC Bind. Buf. 1 - 200 mL	05AR
HTRF Human VHL Bind. Kit - 10K pts Eu Cryptate antibody	02RA
HTRF Human VHL Bind. Kit - 10K pts 6His-VHL complex	02RA
HTRF Human VHL Bind. Kit - 10K pts VHL-Red Ligand	02RA
HTRF Human VHL Bind. Kit - Std	02A

Results :

Incubation : 1h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model : log(inhibitor) vs.response-Variable slope (four parameters)

STD	Conc. (µM)	RATIO
Std 0	0	12113
Std 1	0,0032	11891
Std 2	0,016	11390
Std 3	0,08	9032
Std 4	0,4	5537
Std 5	2	2407
Std 6	10	1361
Std 7	50	1130



Assay characteristics :

Norme provisoire	Found value	Specifications Rev 01
S/B	11	6-12
IC 50	247 nM	150-370 nM

HTRF Human VHL Bind. Kit - 10K pts - Lot n° : 02C

Accepted Rejected

Quality control date : 2024-08-12	Supervisor's initials : MP
-----------------------------------	----------------------------

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.